

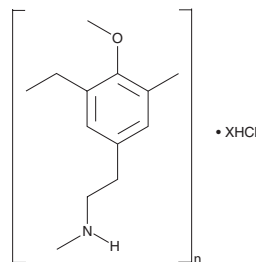
PRODUCT INFORMATION



Compound 48/80 (hydrochloride)

Item No. 22173

MF: $C_{32}H_{45}N_3O_3 \cdot XHCl$ (for trimer)
FW: 519.7
Purity: $\geq 95\%$ (mixture of polymers)
UV/Vis.: λ_{max} : 278 nm
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years
Special Conditions: Store under desiccating conditions



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Compound 48/80 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the compound 48/80 (hydrochloride) in water. The solubility of compound 48/80 (hydrochloride) in water is approximately 100 mM. We do not recommend storing the aqueous solution for more than one day.

Description

Compound 48/80 is a condensation product of N-methyl-*p*-methoxyphenethylamine with formaldehyde that causes histamine degranulation from mast cells.¹ Compound 48/80 also inhibits human platelet aggregation through suppression of calmodulin (CaM) ($IC_{50} = 0.41 \mu g/ml$ for CaM-dependent Ca^{2+} -transporter ATPase activity) and phospholipase C ($IC_{50s} = 2.1$ and $5.0 \mu g/ml$ for cytosolic and particulate phosphatidylinositol-specific activities, respectively).^{2,3}

References

1. Rothschild, A.M. Mechanisms of histamine release by compound 48/80. *Br. J. Pharmac.* **38(1)**, 253-262 (1970).
2. Adamczyk-Engelmann, P. and Gietzen, K. Induction of histamine release and calmodulin antagonism are two distinct properties of compound 48/80. *Cell Calcium* **10(2)**, 93-99 (1989).
3. Bronner, C., Charles, W., Didier, M., *et al.* Compound 48/80 is a potent inhibitor of phospholipase C and a dual modulator of phospholipase A_2 from human platelet. *Biochim. Biophys. Acta* **920(3)**, 301-305 (1987).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/06/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM